

## iFluor™ 350 goat anti-mouse IgG (H+L)

Catalog number: 16440, 16730

Unit size: 200 ug, 1 mg

**Product Details** 

Storage Conditions 2-6°C and kept from light. To extend the shelf-life of this product, add an equal volume of

glycerol to make a final concentration of approximately 50% glycerol and store at -20°C.

Expiration Date 12 months upon receiving

Concentration 1 mg/mL

Formulation PBS, 2 mg/mL BSA

**Unit Details** 

Unit 16440 (200 ug) 16730 (1 mg)

Reconstitution Volume 200 uL ddH<sub>2</sub>O 1 mL ddH<sub>2</sub>O

**Antibody Properties** 

Species Reactivity Mouse

Class Secondary

Clonality Polyclonal

Host Goat

**Chemical Properties** 

Molecular Weight ~150000

**Biological Properties** 

Stabilizer None

Appearance Off-white solid

Preparation Goat anti-mouse IgG (H+L) is produced in goat with pooled total mouse IgG, and affinity purified

with mouse IgG coupled beads. The antibody is conjugated with iFluor™ 350 under optimal

condition.

Application Immunofluorescence (IF), Flow Cytometry (FACS)

Soluble In Water

**Spectral Properties** 

Conjugate iFluor™ 350

Excitation Wavelength 345 nm

## **Applications**

AAT Bioquest's iFluor™ dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 555 and 633 nm). iFluor™ 350 goat anti-mouse IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~345 nm and ~442 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 350 goat anti-mouse IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).